

Version 4e Last updated 30 November 2022

# ab222872

## Human Dipeptidyl Peptidase IV ELISA Kit

For the quantitative measurement of human Dipeptidyl Peptidase IV in plasma, serum and cell culture supernatant samples.

This product is for research use only and is not intended for diagnostic use.

## Table of Contents

|                                     |    |
|-------------------------------------|----|
| 1. Overview                         | 1  |
| 2. Protocol Summary                 | 2  |
| 3. Precautions                      | 3  |
| 4. Storage and Stability            | 3  |
| 5. Limitations                      | 4  |
| 6. Materials Supplied               | 4  |
| 7. Materials Required, Not Supplied | 5  |
| 8. Technical Hints                  | 6  |
| 9. Reagent Preparation              | 7  |
| 10. Standard Preparation            | 9  |
| 11. Sample Preparation              | 11 |
| 12. Assay Procedure                 | 12 |
| 13. Data Analysis                   | 14 |
| 14. Typical Data                    | 15 |
| 15. Typical Sample Values           | 16 |
| 16. Species Reactivity              | 18 |
| 17. Troubleshooting                 | 19 |
| 18. Notes                           | 21 |

# 1. Overview

Human Dipeptidyl Peptidase IV ELISA Kit (ab222872) is designed for the quantitative measurement of Dipeptidyl Peptidase IV (DPP4) in human plasma, serum and cell culture samples.

This assay employs a quantitative sandwich enzyme immunoassay technique that measures human DPP4 in less than 4 hours. A polyclonal antibody specific for human DPP4 has been pre-coated onto a 96-well microplate with removable strips. DPP4 in standards and samples is sandwiched by the immobilized antibody and the biotinylated polyclonal antibody specific for DPP4, which is recognized by a streptavidin-peroxidase conjugate. All unbound material is washed away and a peroxidase enzyme substrate is added. The color development is stopped and the intensity of the color is measured.

Dipeptidyl peptidase 4 (DPP4), also known as adenosine deaminase complexing protein 2 (ADCP-2), T-cell activation antigen CD26 or CD26, belongs to the peptidase S9B family and DPPIV subfamily. It is an intrinsic membrane glycoprotein that comprises 766 amino acids and weighs 110 kDa. Its extracellular cysteine-rich region is necessary for association with collagen, dimer formation and optimal dipeptidyl peptidase activity. Under certain stimuli and chronic inflammation, DPP4 can be released from the membrane, constituting a soluble form by matrix metalloproteases (MMPs). As a serine exopeptidase, DPP4 cleaves and inactivates N-terminal X-proline dipeptides of cytokines, chemokines, and neuropeptides involved in inflammation, immunity, and vascular function. Circulating soluble DPP4 is a novel adipokine that may impair insulin sensitivity in an autocrine and paracrine fashion. DPP4 plays a role in glucose homeostasis through proteolytic inactivation of the intestinal peptide incretins, such as glucose-dependent insulinotropic polypeptide (GIP) and glucagon-like peptide 1 (GLP1), which are major regulators of post-prandial insulin secretion.

## 2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add 50  $\mu$ L standard or sample to appropriate wells. Incubate at room temperature for 2 hours



Wash wells. Add 50  $\mu$ L Biotinylated Antibody to all wells. Incubate at room temperature for 1 hour



Wash wells. Add 50  $\mu$ L Streptavidin-Peroxidase Conjugate to all wells. Incubate at room temperature for 30 minutes



Wash wells. Add 50  $\mu$ L Chromogen Substrate to all wells. Incubate at room temperature for 15 minutes



Add 50  $\mu$ L Stop Solution and read OD at 450 nm

### 3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

### 4. Storage and Stability

Store kit at 4°C immediately upon receipt, apart from the Streptavidin-Peroxidase Conjugate and Biotinylated Antibody, which should be stored at -20°C. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

## 5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

## 6. Materials Supplied

| Item  | Quantity  | Storage Condition |
|---|-----------|-------------------|
| Anti- Human DPP4 coated Microplate (12 x 8 wells) | 96 wells  | +4°C              |
| Human DPP4 Standard                               | 1 vial    | +4°C              |
| Biotinylated Human DPP4                           | 120 µL    | -20°C             |
| 10X Diluent M Concentrate                         | 20 mL     | +4°C              |
| 20X Wash Buffer Concentrate                       | 2 x 30 mL | +4°C              |
| 100X Streptavidin-Peroxidase Conjugate            | 80 µL     | -20°C             |
| Chromogen Substrate                               | 7 mL      | +4°C              |
| Stop Solution                                     | 11 mL     | +4°C              |
| Sealing Tapes                                     | 3         | +4°C              |
| 1X Standard Diluent                               | 2 mL      | +4°C              |

## 7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at 450 nm.
- Deionized water.
- Multi- and single-channel pipettes.
- Tubes for standard dilution.

## 8. Technical Hints

- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Make sure all buffers and solutions are at room temperature before starting the experiment.
- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Make sure you have the right type of plate for your detection method of choice.
- Make sure the heat block/water bath and microplate reader are switched on before starting the experiment.



## 9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells.
- Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

### 9.1 1X Diluent M:

If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the 10X Diluent M Concentrate 1:10 with reagent grade water. Store for up to 30 days at 2-8°C.

### 9.2 Biotinylated Human DPP4 Antibody:

Spin down the antibody briefly and dilute the desired amount of the antibody 1:50 with 1 x Diluent M. Any remaining solution should be frozen at -20°C.

### 9.3 1x Wash Buffer:

If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved. Dilute the 20 x Wash Buffer Concentrate 1:20 with reagent grade water.

### 9.4 1x Streptavidin-Peroxidase Conjugate:

Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 1:100 with 1 x Diluent M. Any remaining solution should be frozen at -20°C.

### 9.5 Anti-Human DPP4 coated Microplate (12 x 8 wells):

Ready to use. Unused microplate wells may be returned to the foil pouch with the desiccant packs and resealed. May be stored for up to 30 days in a vacuum desiccator.

### 9.6 Chromagen Substrate:

Ready to use. Store at +4°C.

### 9.7 Sealing Tapes:

Ready to use. Store at +4°C.

### 9.8 Stop Solution:

Ready to use. Store at +4°C.

## 10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

### 10.1 Reconstitute the DPP4 Stock to generate a 400 ng/mL **Standard Stock**.

- 10.1.1 First consult the DPP4 Standard vial to determine the mass of protein in the vial.
- 10.1.2 Calculate the appropriate volume of 1X Diluent M to add when resuspending the DPP4 Standard vial to produce a 400 ng/mL DPP4 Standard stock by using the following equation:

$C_S$  = Starting mass of DPP4 Standard stock (see vial label) (ng)

$C_F$  = 400 ng/mL DPP4 Standard #1 final required concentration

$V_D$  = Required volume of 1X Diluent M for reconstitution ( $\mu$ L)

Calculate total required volume 1X Diluent M for resuspension:

$$(C_S / C_F) * 1,000 = V_D$$

Example:

**NOTE: This example is for demonstration purposes only. Please remember to check your standard vial for the actual amount of standard provided.**

$C_S$  = 2,40 ng of DPP4 Standard in vial

$C_F$  = 400 ng/mL DPP4 **Standard stock** final concentration

$V_D$  = Required volume of 1X Diluent M for reconstitution

$$(2,40 \text{ ng} / 400 \text{ ng/mL}) * 1,000 = 600 \mu\text{L}$$

- 10.1.3 First briefly centrifuge the DPP4 Standard Vial to collect the contents on the bottom of the tube.
- 10.1.4 Reconstitute the DPP4 Standard vial by adding the appropriate calculated amount VD of 1X Diluent M to the vial to generate the 400 ng/mL DPP4 **Standard stock**. Mix gently and thoroughly.
- 10.2 Allow the reconstituted 400 ng/mL DPP4 **Standard stock** to sit for 10 minutes with gentle agitation prior to making subsequent dilutions
- 10.3 To produce **Standard #1** dilute the standard stock to 200 ng/ml
- 10.4 Label seven tubes **#2 – 8**.
- 10.5 Prepare duplicate or triplicate standard points by serially diluting the **standard #1** solution (200 ng/mL) 1:2 with 1 x Diluent M to produce 100, 50, 25, 12.5, 6.25, 3.125 ng/mL solutions. 1 x Diluent M serves as the zero standard (0 ng/mL). Aliquot standard to limit repeated freezing and thawing. Any remaining solution in the aliquot tube should be frozen at -20°C and used within 3 days. Avoid repeated freeze-thaw cycles.
- 10.6 Add 120 µL of 1X Diluent M to tube **#2 – 8**.
- 10.7 To prepare **Standard #2**, add 120 µL of the **Standard #1** into tube #2 and mix gently.
- 10.8 To prepare **Standard #3**, add 120 µL of the **Standard #2** into tube #3 and mix gently.
- 10.9 Using the table below as a guide, prepare subsequent serial dilutions.

| Standard # | Volume to dilute (µL) | Volume Diluent M (µL) | DPP4 (ng/mL) |
|------------|-----------------------|-----------------------|--------------|
| 1          | Step 10.1             |                       | 200          |
| 2          | 120 µL Standard #1    | 120                   | 100          |
| 3          | 120 µL Standard #2    | 120                   | 50           |
| 4          | 120 µL Standard #3    | 120                   | 25           |
| 5          | 120 µL Standard #4    | 120                   | 12.5         |
| 6          | 120 µL Standard #5    | 120                   | 6.25         |
| 7          | 120 µL Standard #6    | 120                   | 3.125        |
| 8 (Blank)  | N/A                   | 120                   | 0            |

## 11. Sample Preparation

### 11.1 Plasma:

Collect plasma using one-tenth volume of 0.1 M sodium citrate as an anticoagulant. Centrifuge samples at 3000 x g for 10 minutes. Dilute plasma samples 1:4 with 1 x Diluent M and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles. (EDTA or Heparin can also be used as anticoagulant).

### 11.2 Serum:

Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 3000 x g for 10 minutes, and remove serum. Dilute serum samples 1:4 with 1 x Diluent M and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

### 11.3 Cell Culture Supernatants:

Centrifuge cell culture media at 1500 rpm for 10 minutes at 4°C to remove debris. Collect supernatants and assay. If necessary, dilute samples into EIA Diluent; user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.

## 12. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
  - We recommend that you assay all standards, controls and samples in duplicate.
  - Prepare all reagents, working standards, and samples as directed in the previous sections.
- 12.1 Prepare all reagents, working standards, and samples as directed in the previous sections. The assay is performed at room temperature (20-25°C).
  - 12.2 Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccants inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
  - 12.3 Add 50  $\mu$ L of Human DPP4 Standard or sample to each well. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
  - 12.4 Wash five times with 200  $\mu$ L of Wash Buffer manually. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a machine, wash six times with 300  $\mu$ L of Wash Buffer and then invert the plate, decanting the contents; hit 4-5 times on absorbent material to completely remove the liquid.
  - 12.5 Add 50  $\mu$ L of Biotinylated Human DPP4 Antibody to each well and incubate for 1 hour.
  - 12.6 Wash the microplate as described above (Step 12.4).
  - 12.7 Add 50  $\mu$ L of Streptavidin-Peroxidase Conjugate to each well and incubate for 30 minutes.
  - 12.8 Turn on the microplate reader and set up the program in advance.
  - 12.9 Wash the microplate as described above (Step 12.4).
  - 12.10 Add 50  $\mu$ L of Chromogen Substrate to each well and incubate for 15 minutes or until the optimal blue color density develops. Gently tap plate to ensure thorough mixing and break the bubbles in the well with pipette tip
  - 12.11 Add 50  $\mu$ L of Stop Solution to each well. The color will change from blue to yellow.

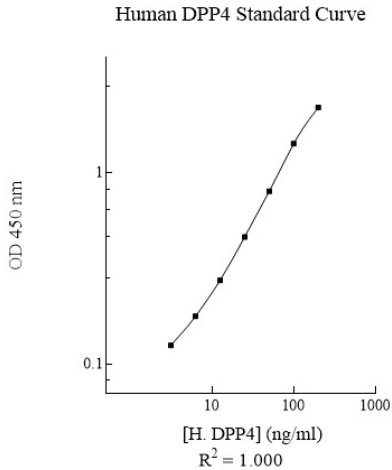
12.12 Read the absorbance on the microplate reader at a wavelength of 450 nm immediately. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. **Δ Note:** Some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

## 13. Data Analysis

- 13.1 Calculate the mean value of the duplicate or triplicate readings for each standard and sample.
- 13.2 To generate a standard curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance (OD) on the y-axis. The best-fit line can be determined by regression analysis using log-log or four-parameter logistic curve-fit.
- 13.3 Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.

## 14. Typical Data

**Typical standard curve** – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.



**Figure 1.** Example of human Dipeptidyl Peptidase IV standard curve.



## 15. Typical Sample Values

### SENSITIVITY –

The minimum detectable dose of DPP4 as calculated by 2SD from the mean of a zero standard was established to be 1.8 ng/mL.

### PRECISION –

Intra-assay precision was determined by testing replicates of three plasma samples in one assay.

Inter-assay precision was determined by testing three plasma samples in twenty assays.

| Sample            | Intra-Assay Precision |     |     | Inter-Assay Precision |      |      |
|-------------------|-----------------------|-----|-----|-----------------------|------|------|
|                   | 1                     | 2   | 3   | 1                     | 2    | 3    |
| n                 | 20                    | 20  | 20  | 20                    | 20   | 20   |
| CV (%) Recovery % | 3.0                   | 3.6 | 3.9 | 8.5                   | 10.1 | 10.5 |
| Average CV (%)    | 3.5                   |     |     | 9.7                   |      |      |

### RECOVERY –

|                      |                  |
|----------------------|------------------|
| Standard Added Value | 12.5 – 100 ng/mL |
| Recovery %           | 91 – 109 %       |
| Average Recovery %   | 96%              |

## Linearity of Dilution

| Average Percentage of Expected Value (%) |        |       |
|--|--------|-------|
| Sample Dilution                          | Plasma | Serum |
| 2  | 91%    | 94%   |
| 4  | 100%   | 101%  |
| 8  | 104%   | 108%  |

## 16. Species Reactivity

This kit recognizes human Dipeptidyl Peptidase IV protein.  
10% FBS in culture media will not affect the assay.

| Species | Cross Reactivity (%) |
|---------|----------------------|
| Canine  | None                 |
| Bovine  | None                 |
| Monkey  | None                 |
| Mouse   | None                 |
| Rat     | None                 |
| Swine   | 5                    |
| Rabbit  | None                 |

### REFERENCE VALUE -

Human plasma and serum samples from healthy adults were tested (n=30). On average, DPP4 level was 0.71 µg/mL.

Please contact our Technical Support team for more information.

## 17. Troubleshooting

| Problem              | Reason                                    | Solution  |
|----------------------|---|---|
| <b>Low Precision</b> | Use of expired components                 | Check the expiration date listed before use.<br>Do not interchange components from different lots.  |
|                      | Improper wash step                        | Check that the correct wash buffer is being used.<br>Check that all wells are dry after aspiration.<br>Check that the microplate washer is dispensing properly.<br>If washing by pipette, check for proper pipetting technique. |
|                      | Splashing of reagents while loading wells | Pipette properly in a controlled and careful manner.  |
|                      | Inconsistent volumes loaded into wells    | Pipette properly in a controlled and careful manner.<br>Check pipette calibration.<br>Check pipette for proper performance.   |
|                      | Insufficient mixing of reagent dilutions  | Thoroughly agitate the lyophilized components after reconstitution.<br>Thoroughly mix dilutions.  |
|                      | Improperly sealed microplate              | Check the microplate pouch for proper sealing.<br>Check that the microplate pouch has no punctures.<br>Check that three desiccants are inside the microplate pouch prior to sealing   |

|  |  |  |
|--|--|--|
| <b>Unexpectedly Low or High Signal Intensity</b> | Microplate was left unattended between steps   | Each step of the procedure should be performed uninterrupted.  |
|  | Omission of step                               | Consult the provided procedure for complete list of steps.   |
|  | Steps performed in incorrect order             | Consult the provided procedure for the correct order.  |
|  | Insufficient amount of reagents added to wells | Check pipette calibration.<br>Check pipette for proper performance.  |
|  | Wash step was skipped                          | Consult the provided procedure for all wash steps.   |
|  | Improper wash buffer                           | Check that the correct wash buffer is being used.  |
|  | Improper reagent preparation                   | Consult reagent preparation section for the correct dilutions of all reagents.   |
|  | Insufficient or prolonged incubation periods   | Consult the provided procedure for correct incubation time.  |
| <b>Deficient Standard Curve fit</b>              | Non-optimal sample dilution                    | Sandwich ELISA: If samples generate OD values higher than the highest standard point (P1), dilute samples further and repeat the assay. User should determine the optimal dilution factor for samples. |
|  | Contamination of reagents.                     | A new tip must be used for each addition of different samples or reagents during the assay procedure.  |
|  | Contents of wells evaporate                    | Verify that the sealing film is firmly in place before placing the assay in the incubator or at room temperature.  |
|  | Improper pipetting                             | Pipette properly in a controlled and careful manner.<br>Check pipette calibration.<br>Check pipette for proper performance.  |
|  | Insufficient mixing of reagent dilutions       | Thoroughly agitate the lyophilized components after reconstitution.<br>Thoroughly mix dilutions.   |

## 18. Notes



## Technical Support

Copyright © 2022 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

**For all technical or commercial enquiries please go to:**

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)